

PRODUCT INFORMATION & MANUAL

Glucose Oxidase Activity Assay Kit (Colorimetric) NBP3-24491

For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Glucose Oxidase Activity Assay Kit (Colorimetric)

Catalog No: NBP3-24491

Method: Colorimetric method

Specification: 96T (Can detect 80 samples without duplication)

Instrument: Microplate reader

Sensitivity: 0.006 U/L

Detection range: 0.006-0.1 U/L

Average intra-assay CV (%): 3

Average inter-assay CV (%): 7

Average recovery rate (%): 101

- ▲ This kit is for research use only.
- ▲ Instructions should be followed strictly, changes of operation may result in unreliable results.
- ▲ Please kindly provide us the lot number (on the outside of the box) of the kit for more efficient service.

General information

▲ Intended use

This kit can be used to measure glucose oxidase (GOD) activity in samples.

▲ Detection principle

Glucose oxidase is an important industrial enzyme in the food industry, widely used in wine, beer, juice, milk powder and other food deoxygenation, flour improvement, food browning prevention and other aspects, in food rapid detection and biosensors are also widely used. Microorganisms are the main source of GOD production with rapid growth and reproduction and wide sources, and the main production strains are aspergillus niger and penicillium.

Glucose oxidase can catalyze the oxidation of glucose to produce hydrogen peroxide. In the presence of peroxidase, hydrogen oxidizes pigment sources to form colored substances. Measure the OD value at 550 nm and glucose oxidase (GOD) activity can be calculated indirectly.

▲ Kit components & storage

| Item | Component | Specification | Storage |
|-----------|------------------------|------------------|----------------------------------|
| Reagent 1 | Buffer Solution | 30 mL × 1 vial | -20°C , 12 months |
| Reagent 2 | Enzyme Reagent | Powder × 2 vials | -20°C , 12 months, shading light |
| Reagent 3 | Chromogenic Agent | 1.6 mL × 2 vials | -20°C , 12 months, shading light |
| Reagent 4 | 1 mmol/L Standard | 1.6 mL × 2 vials | -20°C , 12 months, shading light |
| | Microplate | 96 wells | No requirement |
| | Plate Sealer | 2 pieces | |

Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other.

▲ Materials prepared by users



Instruments

MicroPipettor, Water bath, Centrifuge, Microplate reader (540-560 nm), Incubator, 10 KD Ultrafiltration tube

▲ Safety data

Some of the reagents in the kit contain dangerous substances. It should be avoided to touch the skin and clothing. Wash immediately with plenty of water if touching it carelessly. All the samples and waste material should be treated according to the relevant rules of laboratory's biosafety.

▲ Precautions

Before the experiment, please read the instructions carefully, and wear gloves and work clothes.

Pre-assay preparation

▲ Reagent preparation

- 1. Bring all reagents to room temperature before use.
- 2. Preparation of reagent 2 working solution:

Dissolve a vial of powder with 0.2 mL double distilled water. Prepare fresh solution before use and can be stored at -20°C for 1 week. Avoid to repeated freeze-thawing.

3. The preparation of reaction working solution:

Mix the reagent 1 and reagent 2 working solution at a ratio of 49:1. Prepare the fresh solution before use. Preserve the prepared working solution on the ice for use, and should be used up within 4 h.

▲ Sample preparation

Sample:

Accurately weigh the tissue, add 9 times the volume of normal saline (0.9% NaCl) according to the ratio of Weight (g): Volume (mL) =1:9. Homogenize the tissue sample with homogenizer on ice. Centrifuge the homogenized tissue at 10000 g for 10 min, then take the supernatant and preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant.

▲ Dilution of sample

It is recommended to take 2~3 samples with expected large difference to do pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (0.006–0.1 U/L).

Assay protocol

▲ Plate set up

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|---|---|----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| Α | Α | Α | S1 | S9 | S17 | S25 | S33 | S41 | S49 | S57 | S65 | S73 |
| В | В | В | S2 | S10 | S18 | S26 | S34 | S42 | S50 | S58 | S66 | S74 |
| С | С | С | S3 | S11 | S19 | S27 | S35 | S43 | S51 | S59 | S67 | S75 |
| D | D | D | S4 | S12 | S20 | S28 | S36 | S44 | S52 | S60 | S68 | S76 |
| Е | Е | Е | S5 | S13 | S21 | S29 | S37 | S45 | S53 | S61 | S69 | S77 |
| F | F | F | S6 | S14 | S22 | S30 | S38 | S46 | S54 | S62 | S70 | S78 |
| G | G | G | S7 | S15 | S23 | S31 | S39 | S47 | S55 | S63 | S71 | S79 |
| Н | Н | Н | S8 | S16 | S24 | S32 | S40 | S48 | S56 | S64 | S72 | S80 |

Note: A-H, standard wells; S1-S80, sample wells.

▲ Detailed operation steps

1. The preparation of standard curve

Dilute 1 mmol/L standard with double distilled water to a serial concentration. The recommended dilution gradient is as follows: 0, 0.2, 0.3, 0.4, 0.6, 0.8, 0.9, 1 mmol/L. Reference is as follows:

| Number | Standard concentrations (mmol/L) | 1 mmol/L standard solution (µL) | Double distilled water (µL) |
|--------|----------------------------------|---------------------------------|--------------------------------|
| А | 0 | 0 | 100 |
| В | 0.2 | 20 | 80 |
| С | 0.3 | 30 | 70 |
| D | 0.4 | 40 | 60 |
| Е | 0.6 | 60 | 40 |
| F | 0.8 | 80 | 20 |
| G | 0.9 | 90 | 10 |
| Н | 1 | 100 | 0 |

2. The measurement of samples

- Standard well: Take 10 μL of standard solution with different concentration to the wells.
 - Sample well: Take 10 µL of sample to the wells.
- (2) Add 200 µL of reaction working solution into the each well.
- (3) Add 20 µL of reagent 3 into each well.
- (4) Mix fully for 3 s with microplate reader. Measure the OD value of each well with microplate reader at 550 nm, recorded as A₁.
- (5) Incubate at 37°C for 20 min.
- (6) Measure the OD value of each well with microplate reader at 550 nm, recorded as A_2 , $\Delta A = A_2 A_1$. (Note: There is no change in OD value of standard well, plot the standard curve with the OD value of $A_{2(standard)}$).

▲ Summary operation table

| | Standard well | Sample well |
|--|---------------|-------------|
| Standard with different concentration (µL) | 10 | |
| Sample (µL) | | 10 |
| Reaction working solution (µL) | 200 | 200 |
| Reagent 3 (µL) | 20 | 20 |

Mix fully. Measure the OD value of each well, recorded as A₁.

Incubate at 37°C for 20 min. Mix fully. Measure the OD value of each well, recorded as A_2 , $\Delta A = A_2 - A_1$. (Note: There is no change in OD value of standard well, plot the standard curve with the OD value of $A_{2(standard)}$).

▲ Calculation

Plot the standard curve by using OD value of standard and correspondent concentration as y-axis and x-axis respectively. Create the standard curve with graph software (or EXCEL). The concentration of the sample can be calculated according to the formula based on the OD value of sample. The standard curve is: y=ax+b.

Definition: The enzyme amount of 1 µmol of hydrogen generated by 1 g sample protein per minute at 37°C is defined as 1 unit.

GOD activity (U/gprot) =
$$(\Delta A_{550} - b) \div a \div T \times 1000 \div C_{pr} \times f$$

Note:

y: $OD_{Standard} - OD_{Blank}$ (OD_{Blank} is the OD value when the standard concentration

is 0, standard only detect A₂ value).

x: The concentration of standard.

a: The slope of standard curve.

b: The intercept of standard curve.

 ΔA_{550} : ΔA_{Sample} (A_2 - A_1).

T: Reaction time, 20 min

1000: 1 mmol/L=1000 µmol/L

 C_{pr} : Concentration of protein in sample, gprot/L

f: Dilution factor of sample before test.